

## STAT6 Protein

Full-length human protein expressed in Sf9 cells

**Catalog # S57-30G**

Lot # M012-3

### Product Description

Full-length recombinant human STAT6 was expressed in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC075852](#).

### Gene Aliases

STAT6B, STAT6C, D12S1644, IL-4-STAT

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

### Scientific Background

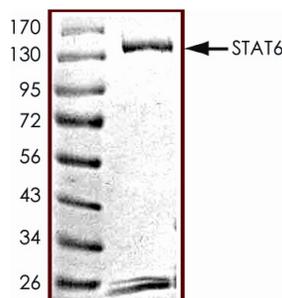
The STAT6 protein is a member of the STAT family of transcription factors that are activated by cytokines and phosphorylated by the receptor associated kinases. STATs form homo- or hetero-dimers that translocate to the cell nucleus where they act as transcription activators (1). STAT6 plays a central role in exerting IL4 mediated biological responses. STAT6 induces the expression of BCL2L1/BCL-XL, which is responsible for the anti-apoptotic activity of IL4. STAT6 mRNA has been detected in various tissues including peripheral blood lymphocytes, colon, intestine, ovary, prostate, thymus, spleen, kidney, liver, lung and placenta. STAT6 is critically involved in Th2 immune response (2).

### References

- Weidinger, S. et al: Association of a STAT 6 haplotype with elevated serum IgE levels in a population based cohort of white adults. *J. Med. Genet.* 41: 658-663, 2004.
- Wang, W. et al: The susceptibility to experimental myasthenia gravis of STAT6<sup>-/-</sup> and STAT4<sup>-/-</sup> BALB/c mice suggests a pathogenic role of Th1 cells. *J. Immun.* 172: 97-103, 2004.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

### Purity



The purity of STAT6 protein was determined to be **>80%** by densitometry.  
Approx. MW **138kDa**.

## STAT6 Protein

Full-length human protein expressed in Sf9 cells

|                     |  |
|---------------------|--|
| Catalog Number      | S57-30G  |
| Specific Lot Number | M012-3   |
| Purity              | >80%   |
| Concentration       | 0.2 µg/µl  |
| Stability           | 1yr at $-70^{\circ}\text{C}$ from date of shipment   |
| Storage & Shipping  | Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice. |